





Quantifoil™ R 1.2/1.3 plus C2 on 200 copper mesh

Cat. No.	Amount
X-101-Cu200C2	100 pcs.

For general laboratory use.

Shipping: shipped at ambient temperature Storage Conditions: store at ambient temperature Shelf Life: n/a

Applications:

Support film for biological samples in Cryo-EM techniques

Description:

QUANTIFOILTM R 1.2/1.3 plus C2 is a 12 nm carbon film with circular holes of 1.2 μ m diameter and an interspace of 1.3 μ m in both dimensions, which has an additional 2 nm continuous carbon layer on top. Applied on 200 copper mesh.

